



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Mouse IL-1 β Kit
Lot Number: K0082964
Expiration Date: 31 DEC 2027
Catalog Number: K152QP-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IL-1 β Antibody	D0082444	2-8°C	31 AUG 2028
Mouse IL-1 β 96-Well Small Spot SECTOR Plate	Z0048271	2-8°C	30 NOV 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080429	2-8°C	31 MAY 2030
Diluent 41	M0410059	$\leq -10^{\circ}\text{C}$	31 DEC 2027
Diluent 45	M0450056	$\leq -10^{\circ}\text{C}$	29 FEB 2028

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-1 β	0.72	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 β	1650	pg/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		15 JUL 2026